

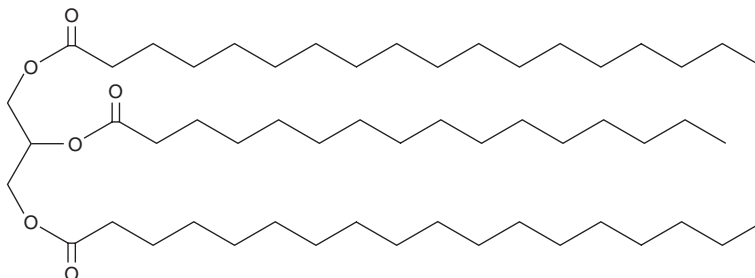
# PRODUCT INFORMATION



## 1,3-Distearoyl-2-Palmitoyl Glycerol

Item No. 27065

**CAS Registry No.:** 2190-24-1  
**Formal Name:** octadecanoic acid, 2-[(1-oxohexadecyl)oxy]-1,3-propanediyl ester  
**Synonyms:** 1,3-Stearin-2-Palmitin, TG(18:0/16:0/18:0)  
**MF:**  $C_{55}H_{106}O_6$   
**FW:** 863.4  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1,3-Distearoyl-2-palmitoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-distearoyl-2-palmitoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Distearoyl-2-palmitoyl glycerol is slightly soluble in chloroform and methanol.

### Description

1,3-Distearoyl-2-palmitoyl glycerol is a triacylglycerol that contains stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-3 positions and palmitic acid (Item No. 10006627) at the *sn*-2 position. It has been found in bovine milk fat.<sup>1</sup>

### Reference

1. Mottram, H.R. and Evershed, R.P. Elucidation of the composition of bovine milk fat triacylglycerols using high-performance liquid chromatography-atmospheric pressure chemical ionisation mass spectrometry. *J. Chromatogr. A*. **926**(2), 239-253 (2001).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)